

# pdUp

## Nucleoside bisphosphate suitable for the characterisation of ribonucleases and nucleotidases

pdUp is a nucleoside bisphosphate that is suitable for characterising ribonucleases, nucleotidases and other nucleotide-binding enzymes. pdUp is accepted as a substrate by T4 RNA ligase for incorporation into oligodeoxyribonucleotides.

*BLG-D262-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

### Ordering Information [Order Online »](#)

BLG-D262-05	5µmol
BLG-D262-25	5x5µmol

### Manuals, SDS & CofA [View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name 2'-Deoxyuridine-3',5'- O- bisphosphate . sodium salt

Appearance Solid

CAS 29774-37-6

Couple Type Modified nucleotides

Formula  $C_9H_{14}N_2O_{11}P_2$  (free acid)

MW 388.2 (free acid)

Purity > 95 % HPLC

Quantity 5 µmol are equivalent to ~1.9 mg

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)